

Ro 41-5253

Retinoic acid receptor ligand

A selective RAR α antagonist (concentration required to inhibit 50% of specific retinoic acid binding: RAR α , 60 nM; RAR β , 2400 nM; RAR γ , 3300 nM). Counteracts the effects of retinoic acid on HL-60 cell differentiation, B-lymphocyte polyclonal activation and in a RAR α transactivation assay. Ro 41-5253 suppresses retinoic acid-induced connexin 43 protein expression in 10T1/2 cells.

Citations: 9

[View Online »](#)

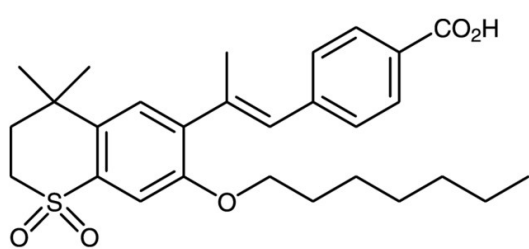
Ordering Information

[Order Online »](#)

BML-GR110-0005	5mg
BML-GR110-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 144092-31-9

Couple Target Retinoic acid receptor

Couple Type Ligand

Formula $C_{28}H_{36}O_5S$

MW 484.7

Purity ≥98% (TLC)

Solubility Soluble in DMSO (50mg/ml) or 100% ethanol (15mg/ml, warm).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)